

PRODUCT INFORMATION



M-CSF- γ (human, recombinant)

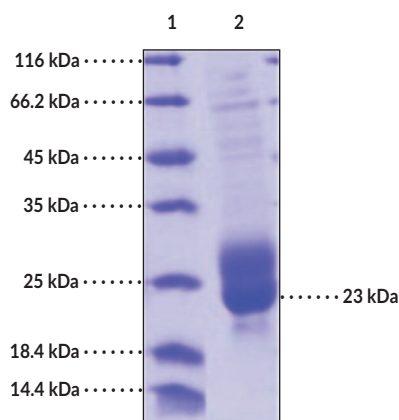
Item No. 32063

Overview and Properties

Synonym:	CSF-1
Source:	Active recombinant human M-CSF- γ expressed in HEK293 cells
Amino Acids:	37-190
Uniprot No.:	P09603-2
Molecular Weight:	18.4 kDa
Storage:	-80°C (as supplied)
Stability:	≥ 1 year
Purity:	$\geq 85\%$ estimated by SDS-PAGE
Supplied in:	Lyophilized from sterile PBS, pH 7.4
Endotoxin Testing:	< 1.0 EU/ μ g, determined by the LAL endotoxin assay
Bioactivity:	See figures for details

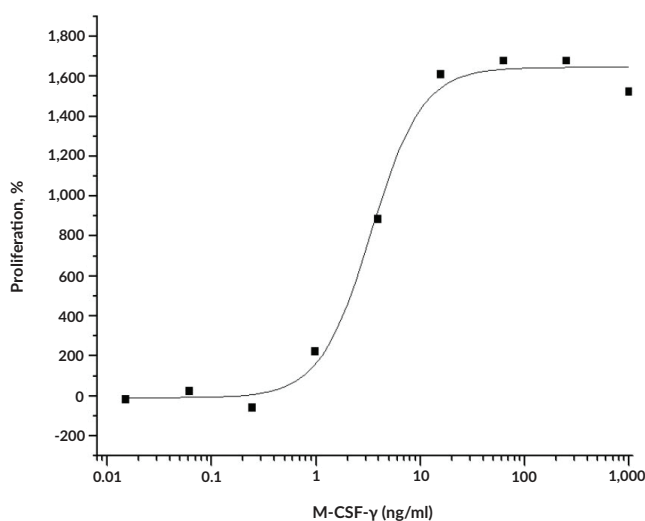
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Images



Lane 1: MW Markers
Lane 2: M-CSF- γ

SDS-PAGE Analysis of M-CSF- γ . This protein has a calculated molecular weight of 18.4 kDa. It has an apparent molecular weight of approximately 23 kDa by SDS-PAGE under reducing conditions due to apparent post-translational modifications.



Cell Proliferation Assay Using M-NFS-60 Mouse Myelogenous Leukemia Lymphoblast Cells. The ED₅₀ value for this effect is typically 1.6-6.3 ng/ml.

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 06/04/2020

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Description

Macrophage colony-stimulating factor (M-CSF) is a glycoprotein encoded by the *CSF1* gene in humans that promotes the differentiation, proliferation, and function of mononuclear phagocytes, including macrophages, osteoclasts, and dendritic cells.^{1,2} Alternative splicing of *CSF1* pre-mRNA produces one full-length long isoform (M-CSF- β), an intermediate-length isoform (M-CSF- γ), and a short-length isoform (M-CSF- α) that share sequence homology in the 150-amino acid receptor binding region that is required for the biological activity of M-CSF.^{3,4} M-CSF exists as a disulfide-linked homodimer where each monomer contains four α -helices, an antiparallel β -sheet, and numerous glycosylation sites.¹ M-CSF is constitutively produced by many cell types, including stromal cells, osteoclasts, fibroblasts, and macrophages, and is localized to the cell surface where it can be proteolytically cleaved to yield a secreted form.^{5,6} Binding of M-CSF to the M-CSF receptor, which is expressed by monocytes, macrophages, osteoclasts, and dendritic cells, promotes cell differentiation, proliferation, and survival of mononuclear phagocytes and regulates bone resorption by osteoclasts.^{5,7} Mice homozygous for *Csf1^{op}*, an inactivating mutation, exhibit defects in fertility and neural development and develop osteopetrosis, a condition characterized by increased bone density.⁸ Neutralization of M-CSF with a monoclonal antibody decreases joint swelling and distortion in a mouse model of collagen-induced arthritis.⁹ Serum M-CSF levels are increased in patients with colorectal, pancreatic, prostate, or head and neck cancer.⁵ M-CSF has been used to generate bone marrow-derived macrophages with an anti-inflammatory phenotype *in vitro*.² Cayman's M-CSF- γ (human, recombinant) protein can be used for cell-based assay applications. This protein consists of 158 amino acids, has a calculated molecular weight of 18.4 kDa, and a predicted N-terminus of Glu37 after signal peptide cleavage. By SDS-PAGE, under reducing conditions, the molecular mass of the protein is approximately 23 kDa due to apparent post-translational modifications.

References

1. Stanley, E.R., Berg, K.L., Einstein, D.B., *et al.* Biology and action of colony-stimulating factor-1. *Mol. Reprod. Dev.* **46(1)**, 4-10 (1997).
2. Ushach, I. and Zlotnik, A. Biological role of granulocyte macrophage colony-stimulating factor (GM-CSF) and macrophage colony-stimulating factor (M-CSF) on cells of the myeloid lineage. *J. Leukoc. Biol.* **100(3)**, 481-489 (2016).
3. Cerretti, D.P., Wignall, J., Anderson, D., *et al.* Human macrophage-colony stimulating factor: Alternative RNA and protein processing from a single gene. *Mol. Immunol.* **25(8)**, 761-770 (1988).
4. Fitzgerald, K.A., Callard, R.E., O'Neill, L.A.J., *et al.* M-CSF. *The Cytokine Factsbook and Webfacts*. Fitzgerald, K.A., editor, 2nd edition, *Academic Press* (2001).
5. Chockalingam, S. and Ghosh, S.S. Macrophage colony-stimulating factor and cancer: A review. *Tumour Biol.* **35(11)**, 10635-10644 (2014).
6. Hamilton, J.A. Colony-stimulating factors in inflammation and autoimmunity. *Nat. Rev. Immunol.* **8(7)**, 533-544 (2008).
7. Stanley, E.R. and Chitu, V. CSF-1 receptor signaling in myeloid cells. *Cold Spring Harb. Mol. Case Stud.* **6(6)**, a021857 (2014).
8. Dai, X.-M., Ryan, G.R., Hapel, A.J., *et al.* Targeted disruption of the mouse colony-stimulating factor 1 receptor gene results in osteopetrosis, mononuclear phagocyte deficiency, increased primitive progenitor cell frequencies, and reproductive defects. *Blood* **99(1)**, 111-120 (2002).
9. Campbell, I.K., Rich, M.J., Bischof, R.J., *et al.* The colony-stimulating factors and collagen-induced arthritis: Exacerbation of disease by M-CSF and G-CSF and requirement for endogenous M-CSF. *J. Leukoc. Biol.* **68(1)**, 144-150 (2000).

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM